

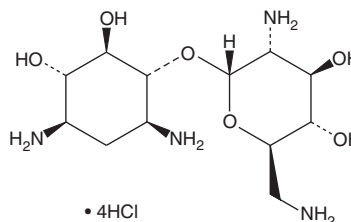
PRODUCT INFORMATION



Neamine (hydrochloride)

Item No. 21752

CAS Registry No.: 15446-43-2
Formal Name: 2-deoxy-4-O-(2,6-diamino-2,6-dideoxy- α -D-glucopyranosyl)-D-streptamine, tetrahydrochloride
Synonym: Neomycin A
MF: $C_{12}H_{26}N_4O_6 \cdot 4HCl$
FW: 468.2
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Neamine (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of neamine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of neamine (hydrochloride) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Neamine is a natural degradation product of neomycin (Item No. 14287) and is a common core molecule used in the synthesis of aminoglycoside antibiotics.¹ Formulations containing neamine are used as antibiotic agents.

Reference

1. Greenberg, W.A., Priestley, E.S., Sears, P.S., *et al.* Design and synthesis of new aminoglycoside antibiotics containing neamine as an optimal core structure: Correlation of antibiotic activity with *in vitro* inhibition of translation. *J. Am. Chem. Soc.* **121(28)**, 6527-6541 (1999).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/22/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM